

A Comparison of Two Commercially Available PCR Assays for the Detection of *Vibrio* from Seafood Samples

Annette Hughes¹, David Crabtree¹, Laura Vaahtoranta², Hanna Lehmusto², Matthew Hahs³. Thermo Fisher Scientific, ¹Basingstoke, UK, ²Vantaa, Finland, ³Lenexa, USA.

INTRODUCTION

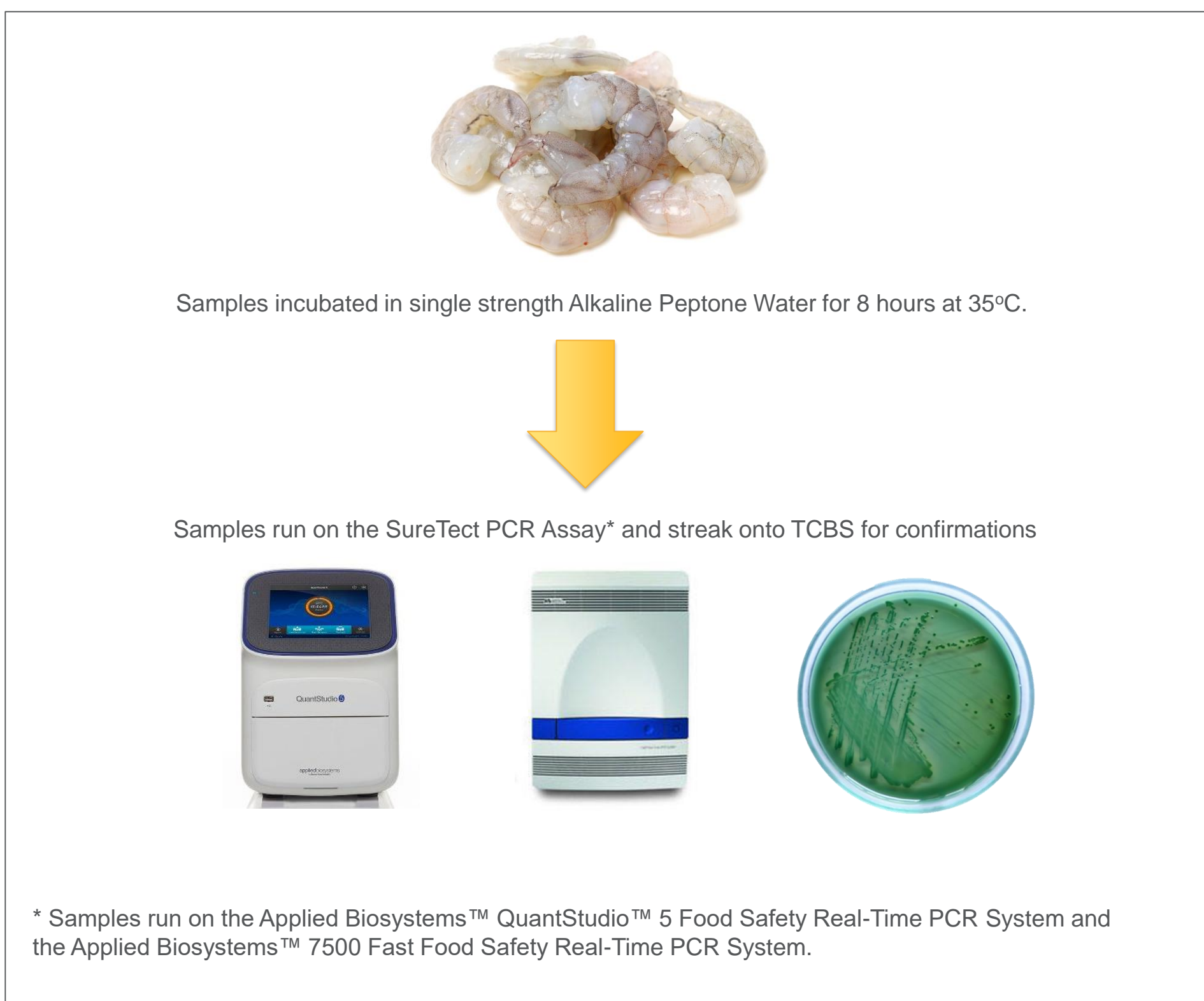
Bivalve mollusks and crustaceans are natural reservoirs of *Vibrio* species and, if eaten raw or under-cooked, carry the risk of causing foodborne disease. A reliable detection method is a valuable tool to ensure that seafood is safe to eat.

This study compares the performance of the Thermo Scientific™ SureTect™ *Vibrio cholerae*, *V. parahaemolyticus* and *V. vulnificus* PCR Assay (candidate method) for detection of target strains from mollusks and crustaceans to the Hygiena™ BAX® System Real-Time PCR Assay for *Vibrio*.

MATERIALS AND METHODS

An unpaired study was conducted testing eighty-eight mollusk and crustacean samples which were artificially contaminated with 2 CFU/25g of either *V. cholerae*, *V. parahaemolyticus* and *V. vulnificus*. Candidate method samples were tested according to the kit workflow (Figure 1) and BAX samples were tested according to the manufacturer's protocol.

Figure 1: Candidate method



RESULTS

Figure 2: Candidate and alternative workflow comparison



Table 1: Method agreement results between the candidate and alternative method

Method agreement	<i>Vibrio cholerae</i>	<i>Vibrio parahaemolyticus</i>	<i>Vibrio vulnificus</i>
Positive agreement	22	27	22
Negative agreement	51	16	45
Positive deviation	9	36	12
Negative deviation	6	9	9

Key:

Positive Agreement = Candidate Method Positive, Alternative Method Positive
 Negative Agreement = Candidate Method Negative, Alternative Method Negative

Positive Deviation = Candidate Method Positive, Alternative Method Negative
 Negative Deviation = Candidate Method Negative, Alternative Method Positive

The results show that the candidate method had superior performance when detecting the different *Vibrio* targets compared to the alternative method, achieving a total of 57 positive deviations on both the QuantStudio 5 and the 7500 Fast (Table 1). The candidate method achieved a time to result of less than 10 hours, compared to 20-22 hours for the BAX method (Figure 2).

CONCLUSIONS

Faster Time to Result

The SureTect workflow delivered more accurate results in a shorter timeframe
 SureTect workflow: <10 hours
 BAX workflow: 20-22 hours

Improved Sensitivity

Improved performance in detection of *V. cholerae*, *V. parahaemolyticus* and *V. vulnificus* from mollusks and crustaceans compared to the BAX PCR Assay.

Simpler workflow

With single enrichment, pre-filled lysis reagents and simple lysis the SureTect workflow delivers more accurate results with a simpler workflow compared to the BAX workflow.

TRADEMARKS/LICENSING

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. BAX® is the property of Hygiena International Ltd. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

LT2586A
 October 2020